

# PRODUCT INFORMATION

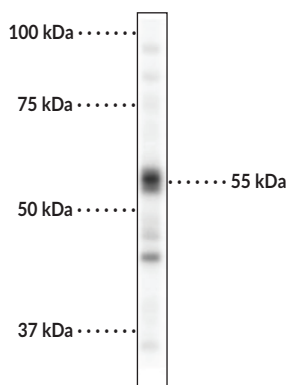


## GABA<sub>A</sub> Receptor $\alpha_5$ Subunit Polyclonal Antibody Item No. 29271

### Overview and Properties

<b>Contents:</b>	This vial contains 100 $\mu$ l of affinity-purified rabbit polyclonal antibody.
<b>Synonyms:</b>	$\gamma$ -Aminobutyric Acid Receptor Subunit $\alpha_5$ , GABA <sub>A</sub> Receptor Subunit $\alpha_5$ , GABRA5
<b>Immunogen:</b>	Fusion protein from the cytoplasmic loop of the $\alpha_5$ subunit of the rat GABA <sub>A</sub> receptor
<b>Molecular Weight:</b>	~55 kDa
<b>Species Reactivity:</b>	(+) Mouse, rat
<b>Form:</b>	Liquid
<b>Storage:</b>	-20°C (as supplied)
<b>Stability:</b>	$\geq$ 1 year
<b>Storage Buffer:</b>	10 mM HEPES, pH 7.5, with 150 mM sodium chloride, 100 $\mu$ g/ml BSA, and 50% glycerol
<b>Host:</b>	Rabbit
<b>Applications:</b>	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution for WB is 1:500. IHC and other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

### Image



WB of mouse whole brain showing specific immunolabeling of the ~55 kDa  $\alpha_5$ -subunit of the GABA<sub>A</sub> receptor.

**WARNING**  
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

**SAFETY DATA**  
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

**WARRANTY AND LIMITATION OF REMEDY**  
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 06/22/2020

**CAYMAN CHEMICAL**  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM

# PRODUCT INFORMATION



## Description

---

GABA<sub>A</sub> receptors are ligand-gated chloride channels that mediate the effects of the inhibitory neurotransmitter GABA in the CNS.<sup>1,2</sup> They are postsynaptic heteropentameric receptors that contain protein subunits from the following isoforms:  $\alpha_{1-6}$ ,  $\beta_{1-4}$ ,  $\gamma_{1-3}$ ,  $\delta$ ,  $\epsilon$ ,  $\pi$ ,  $\theta$ , and  $\rho_{1-3}$ , arranged around a central pore. Phasic inhibitory synaptic transmission is regulated by  $\alpha_1\beta_2\gamma_2$  subunit-containing GABA<sub>A</sub> receptors, the major isoform found in the brain.<sup>2,3</sup> The  $\alpha$  subunit of GABA<sub>A</sub> receptors interfaces with a  $\beta$  subunit to form the GABA binding site that initiates GABA-induced action potentials and forms the benzodiazepine binding site with the  $\gamma$  subunit. Approximately 5% of all GABA<sub>A</sub> receptors contain  $\alpha_5$  subunits and up to 25% of these receptors are expressed in the hippocampus, a region of the brain associated with learning and memory.<sup>4</sup> Deletions in *GABRA5*, which encodes the  $\alpha_5$  subunit isoform, have been found on chromosome 15 in patients with Angelman syndrome who typically develop autism spectrum disorder (ASD). Levels of  $\alpha_5$  subunit-containing GABA<sub>A</sub> receptors are decreased in postmortem hippocampal samples from patients with advanced Alzheimer's disease, Down syndrome, or fetal alcohol syndrome. Cayman's GABA<sub>A</sub> Receptor  $\alpha_5$  Subunit Polyclonal Antibody can be used for immunohistochemistry (IHC) and Western blot (WB) applications. The antibody recognizes the GABA<sub>A</sub> receptor  $\alpha_5$  subunit at approximately 55 kDa from mouse and rat samples.

## References

---

1. Crestani, F. and Rudolph, U. Behavioral functions of GABA<sub>A</sub> receptor subtypes - the Zurich experience. *Adv. Pharmacol.* **72**, 37-51 (2015).
2. Hirose, S. Mutant GABA<sub>A</sub> receptor subunits in genetic (idiopathic) epilepsy. *Prog. Brain Res.* **213**, 55-85 (2014).
3. Wongsamitkul, N., Maldifassi, M.C., Simeone, X., et al.  $\alpha$  subunits in GABA<sub>A</sub> receptors are dispensable for GABA and diazepam action. *Sci. Rep.* **7(1)**, 15498 (2017).
4. Mohamad, F.H., and Has, A.T.C. The  $\alpha_5$ -containing GABA<sub>A</sub> receptors-a brief summary. *J. Mol. Neurosci.* **67(2)**, 343-351 (2019).

CAYMAN CHEMICAL  
1180 EAST ELLSWORTH RD  
ANN ARBOR, MI 48108 · USA  
PHONE: [800] 364-9897  
[734] 971-3335  
FAX: [734] 971-3640  
CUSTSERV@CAYMANCHEM.COM  
WWW.CAYMANCHEM.COM